

<b>Product Number:</b>	101735
<b>Lot Number:</b>	501718
<b>Expiration:</b>	22-Jan-2020

Component	Item Number	Lot Number
SBG Concentrate	103397	907013
SBG Dilution Buffer	100376	902217

SBG Concentrate Concentration (nM)	50
------------------------------------	----

Review/Approval Patrizia Stadler Associate Director, QC *p-Stadler* 18 Mar 2019

Name Title Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.